

Data Sheet

LILRB4, Avi-His-Tag, Biotin-Labeled

Human, Recombinant, C-terminal Avi-His-Tag
Catalog #: 100336
Lot#: 190403 **Conc.:** 1.0 mg/ml

Formulated in: 8 mM phosphate, pH 7.4,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Stability: At least 6 months at -80°C . *Avoid
freeze/thaw cycles.*

References:

1. Brown, D., *et al.*, *BMC Immunology*.
2009 Oct 2009; **10(56)**: 1924-1937.
2. Cheng, H., *et al.*, *J Biol Chem*. May
2011; **286(20)**: 18013-25.

Description: Human LILRB4, also known as leukocyte immunoglobulin like receptor B4, and CD85K, GenBank Accession No. NM_001278426, a.a 17-257, with C-terminal Avi-His-Tag, expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW = 31 kDa. This protein runs at a higher molecular weight due to glycosylation.

Quality Assurance

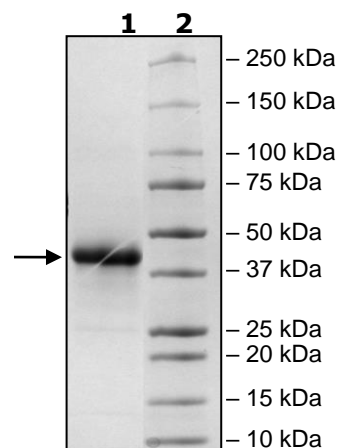
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg LILRB4

Lane 2:
Protein Marker

MW: 31 kDa
+glycans

Purity: $\geq 90\%$



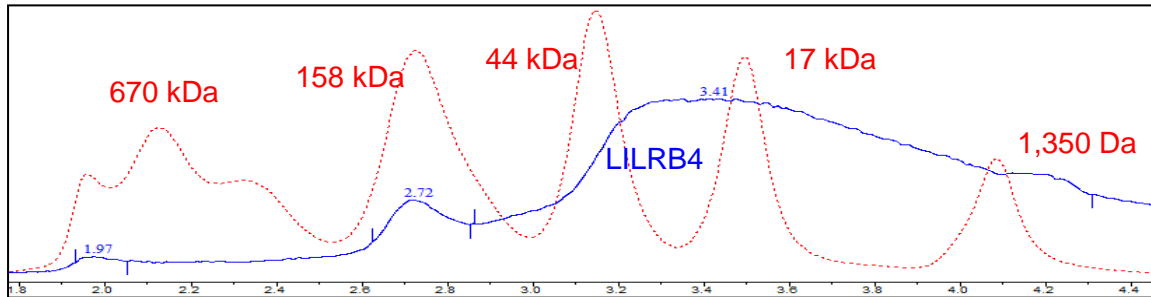
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration:



<10% Aggregation

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com